

4th Joint PhD Student Workshop
January 23/25-2006
Astoria Park Hotel
Riva del Garda (TN)

Monday, 23 January 2006

11:00 - 12:30 Arrival, registration, poster positioning

13:00 - 14:30 Welcome buffet

14:50 - 15:00 Opening Remarks
Andrea Musacchio, IFOM-IEO Campus

15:00 - 16:20 **Cell & Molecular biology 1**

15:00-15:20	<u>Boggio</u>	(IFOM-IEO)	GAM1 AND THE SUMO PATHWAY
15:20-15:40	<u>Crotti</u>	(HRS-DIBIT)	UNDERSTANDING THE ROLE OF STAT5 AND STAT5ΔC IN HIV INFECTION/REPLICATION
15:40-16:00	<u>Ciferri</u>	(IFOM-IEO)	MOLECULAR ARCHITECTURE OF THE HUMAN NDC80-HEC1 COMPLEX, A CRITICAL CONSTITUENT OF THE OUTER KINETOCHORE
16:00-16:20	<u>Marturano</u>	(HRS-DIBIT)	TUMOR ANTIGEN PROCESSING AND PRESENTATION ON MHC CLASS II MOLECULES FOR CD4 ⁺ T CELL RECOGNITION IN HEALTH AND DISEASE

16:20 - 16:50 Coffee break

16:50 - 17:30 **Cell & Molecular biology 2**

16:50-17:10	<u>Sironi</u>	(HRS-DIBIT)	FUNCTIONAL AND BIOCHEMICAL CHARACTERIZATION OF EFA6, A KEY PLAYER IN CYTOSKELETAL REMODELLING
17:10-17:30	<u>Madsen</u>	(IFOM-IEO)	COMPLETE ALANINE-SCAN OF HUMAN UPAR

17:30 - 18:10 **Linda M. Hendershot**

Department of Genetics and Tumor Cell Biology, St. Jude Children's Research Hospital, Memphis, TN, USA

18:10 - 19:00 Round Table

20:00 - 21:30 Dinner

21:30 - 23:00 Poster Session 1

Tuesday, 24 January 2006

9:00 - 11:30 **Cell & Molecular biology 3 (Mix Session)**

- 09:00-09:20 Lombardo (HSR-TIGET) DEVELOPMENT OF NOVEL LENTIVIRAL VECTORS FOR GENE CORRECTION OF INHERITED MUTATIONS
- 09:20-09:40 Biressi (HSR-DIBIT) THE MOLECULAR BASIS OF SKELETAL MYOBLAST DIVERSIFICATION. GENE EXPRESSION ANALYSIS OF PURIFIED EMBRYONIC AND FETAL MYOBLASTS.
- 09:40-10:00 Maiolica (IFOM-IEO) PROTEIN-PROTEIN INTERACTION BY MASS SPECTROMETRY
- 10:00-10:20 Veglianesse (MNI) CHARACTERIZATION OF THE AMPA RECEPTORS SUBUNITS AND THEIR ASSOCIATED TRAFFICKING PROTEINS (ABP, NSF AND PICK-1) IN A MOUSE MODEL OF AMYOTROPHIC LATERAL SCLEROSIS.

10:20 - 10:50 Coffee Break

- 10:50-11:10 Carrassa (MNI) MOLECULAR MECHANISMS REGULATING THE G2 CHECKPOINT FOCUSING ON CHK1
- 11:10-11:30 Magistroni (MNI) BIM PROMOTER IS HIGHLY METHYLATED IN MALIGNANT LYMPHOID CELL LINES, LEADING TO DOWNREGULATION OF BIM EXPRESSION AND PROTECTION FROM APOPTOSIS

11:30 - 13:00 Poster Session 2

13:00 - 14:30 Lunch

14:30 - 16:30 **Immunology Session**

- 14:30-14:50 Baruah (HSR-DIBIT) INNATE OPSONINS: MODULATING PERCEPTION OF CELL DEATH AND INFLAMMATION
- 14:50-15:10 Huber (INT) HUMAN CRC CELLS INDUCE T-CELL DEATH THROUGH RELEASE OF PROAPOPTOTIC MICROVESICLES: ROLE IN IMMUNE ESCAPE
- 15:10-15:30 Cassani (HSR-TIGET) INVESTIGATING THE MOLECULAR MECHANISMS OF IMMUNE DEFICIENCY IN ADA-SCID: IMPLICATIONS FOR GENE THERAPY APPROACH
- 15:30-15:50 Butti (HSR) IL-4 GENE DELIVERY INDUCES CLINICAL AND FUNCTIONAL RECOVERY FROM EXPERIMENTAL AUTOIMMUNE ENCEPHALOMYELITIS, AN ANIMAL MODEL OF MULTIPLE SCLEROSIS.
- 15:50-16:10 Martinoli (IFOM-IEO) DEVELOPMENT AND FUNCTION OF MUCOSAL IGAS INDUCED FOLLOWING ORAL INFECTION WITH *SALMONELLA TYPHIMURIUM*
- 16:10-16:30 Pittoni (HSR) THYMIC SELECTION OF NKT_i CELLS

16:30 - 17:00 Coffee Break

17:00 - 18:20 **Cancer Biology 1**

- 17:00-17:20 Di Micco (IFOM-IEO) RAS-INDUCED SENESCENCE OCCURS VIA A DNA-DAMAGE RESPONSE THAT IS DEPENDENT ON REACTIVE OXYGEN SPECIES-MEDIATED HYPERPROLIFERATION AND ON DNA REPLICATION

17:20-17:40	<u>Bonetti</u>	(IFOM-IEO)	NUCLEOPHOSMIN IS REQUIRED FOR DNA INTEGRITY AND p19ARF PROTEIN STABILITY
17:40-18:00	<u>Polato</u>	(MNI)	CHARACTERIZATION OF A GENE IMPORTANT FOR P53 DEPENDENT, STRESS INDUCED APOPTOSIS
18:00-18:20	<u>Brambilla</u>	(INT)	INTERFERENCE WITH MECHANISMS OF TELOMERE MAINTENANCE BY TARGETING hPOT1 IN HUMAN CANCER CELLS.
20:00	Social Dinner		

Wednesday, 25 January 2006

- 9:00 - 9:40** **Anton Berns**
Netherlands Cancer Institute, Division of Molecular Genetics and Centre of Biomedical Genetics, Amsterdam, The Netherlands.
- 9:40 - 10:30** Round Table
- 10:30 - 11:50** **Cancer Biology 2**
- 10:30-10:50 Cicalese (IFOM-IEO) BIOLOGICAL CHARACTERIZATION OF MOUSE BREAST STEM CELLS FROM NORMAL AND NEOPLASTIC TISSUES
- 10:50-11:10 Cambiaghi (IFOM-IEO) PRELEUKEMIC PHASE ANALYSIS TO UNDERSTAND THE BIOLOGICAL CONTRIBUTION OF PML-RAR α TO THE LEUKEMOGENIC PROCESS
- 11:10- 11:30 De Franco (IFOM-IEO) ROLE OF QUIESCIENT STEM CELLS IN THE MAINTENANCE OF LEUKAEMIA CELL POPULATIONS
- 11:30- 11:50 Ortensi (IFOM-IEO) RAI (SHC-C) REGULATES NEURONAL STEM CELL FUNCTION
- 11:50 - 12:20** Coffee Break
- 12:20 - 13:00** **Cancer Biology 3**
- 12:20 -12:40 Sangaletti (INT) STROMAL-DERIVED SPARC REGULATES STROMA CELLS TUMOR CELLS INTERACTION TOWARD METASTASIS.
- 12:40 -13:00 Maddaluno (IFOM-IEO) L1CAM: A NEW REGULATOR OF DENDRITIC CELL TRAFFICKING
- 13:00 - 14:30** Lunch
- 14:30 - 16:00** Poster Session 3
- 16:00 - 17:30** Science& Society Session
- 16:00 - 16:40** **Lorenzo Tomatis, M.D., Writer**
- 16:40 - 17:30** Round Table
- 17:30 - 17:40** Closing Remarks
Domenico Delia, Istituto Nazionale Tumori